

Fa2N-4 Lot No. 2310144 **Immortalized Human Hepatocytes**

Produced under license from Corning Incorporated

IFH15 – Cryopreserved cells	1.5 mL vial
IFHP06 – Plated cells	6-well plate
IFHP12 – Plated cells	12-well plate
IFHP24 – Plated cells	24-well plates
IFHP48 – Plated cells	48-well plates
IFHP96 – Plated cells	96-well plates

Enzyme	Prototypical Inducer	Marker Substrate Reaction	Fold Induction
CYP1A2	Omeprazole (60 μ M)	Phenacetin O-dealkylation	11.45
CYP2B6	Phenobarbital (750 μ M)	Bupropion hydroxylation	1.08
CYP2C9	Rifampin (40 μ M)	Diclofenac 4'-hydroxylation	3.31
CYP3A4/5	Rifampin (40 μ M)	Midazolam 1'-hydroxylation	1.98
CYP3A4/5	Rifampin (40 μ M)	Atorvastatin hydroxylation	2.82

Fa2N-4 immortalized hepatocytes were plated in 24-well plates at 0.33×10^6 cells per well on Day 0 and cultured in MFE Support Medium F containing Supplement A on days 1-3. Cells were then treated with inducers (in MFE Support Medium F containing Supplement A) once daily for 3 consecutive days. Substrate incubations were performed for 15 minutes to 6 hours. CYP activity was measured under conditions where the metabolite formation was proportional to incubation time. Samples were analyzed by LC/MS/MS. All substrate concentrations were at 100 μ M, except Atorvastatin which was at 250 μ M.

Cell Line Information

The Fa2N-4 cell line was prepared by immortalizing hepatocytes from a 12 year-old Caucasian female donor with the SV40 large T antigen. The donor tested negative for CMV, HIV, HBV, and HCV.



Store in liquid nitrogen, vapor phase

CAUTION: This sample should be considered as a potential biohazard and universal precautions should be followed. Intended for *in vitro* use only.

These data were generated by and are the property of XenoTech. These data are not to be reproduced, published or distributed without the express written consent of XenoTech.

This data sheet serves as a Certificate of Analysis and has been approved by **Stephanie Helmstetter, Assistant Director.**

Signature and Date: Stephanie Helmstetter 22 June 2023